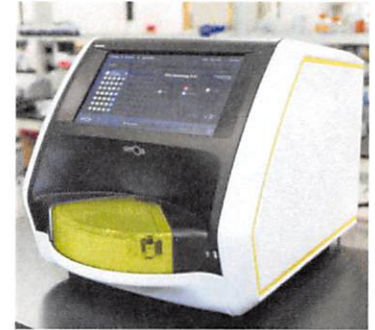


Product Number:	103400
Lot Number:	501964
Expiration:	15-Apr-2020
Platform(s):	HD-1/HD-X and SR-X



Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

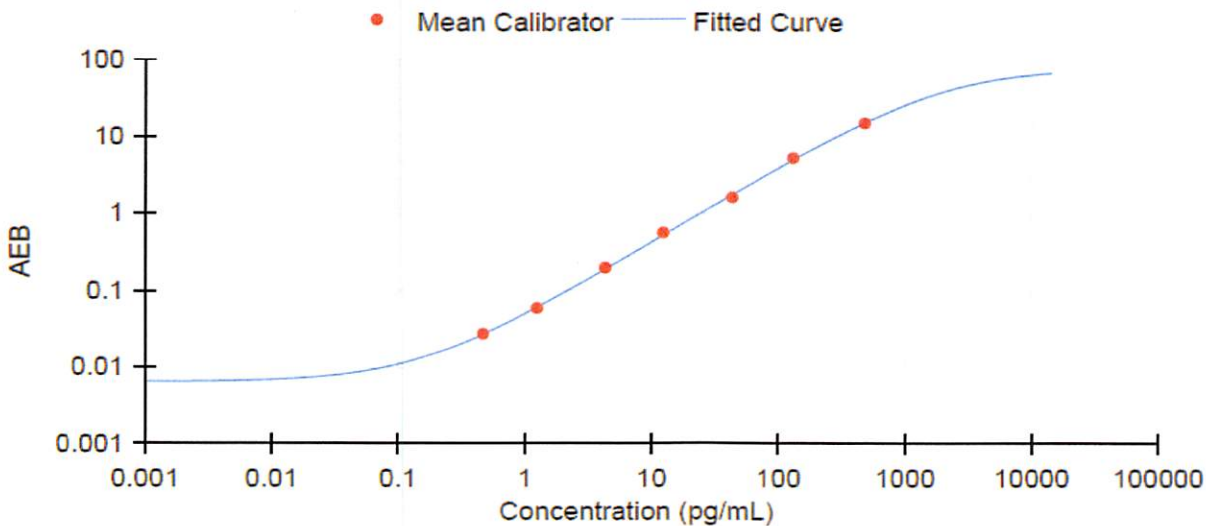
Component	Part Number	Lot Number
Bead Reagent	102246	911413
Detector Reagent	102248	910508
SBG Reagent	102250	910509
Sample Diluent	103399	910505
Calibrators	102255	918901
Controls	102256-102257	918902
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

Release Materials			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918902	3.20	2.30-4.01
Control 2	918902	126	102-154
Panel 1	823301	4.73	4.08-6.68
Panel 2	823302	6.59	4.90-7.61
Panel 3	823601	82.9	59.3-101
Panel 4	823602	334	236-402

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.466
C	1.24
D	4.27
E	12.3
F	42.8
G	130
H	475

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p. Stadler</i> 15 Aug 2019
	Name	Title	Signature/Date